

Product Number:	102732
Lot Number:	501323
Expiration:	3-Jan-2019
Platform(s):	HD-1 & SR-X



Component	Part Number	Lot Number
Bead Reagent	102722	819303
Detector Reagent	102724	818712
SBG Reagent	102726	818709
Sample Diluent	102728	818415
Calibrator Diluent	102731	818711
Calibrator Concentrate	100658	73217
RGP Reagent	103159	N/A ¹

Data below represents results generated on the Simoa™ HD-1 Analyzer.

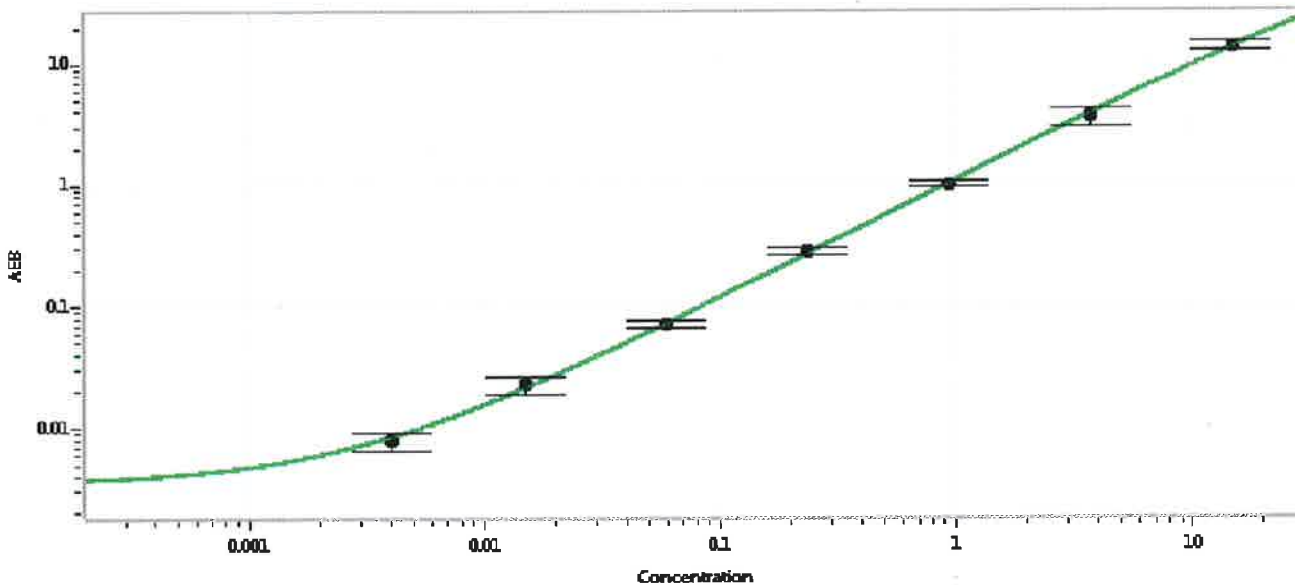
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-061217-01	0.169	0.134-0.289
Control 2	MF-061217-01	3.81	2.44-6.35
Panel 1	MF-061317-03	0.049	0.037-0.088
Panel 2	MF-061317-04	0.280	0.191-0.490
Panel 3	MF-061317-05	4.94	3.18-7.89

* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibrator Concentrate Value Assignment (pg/mL)	194
-------------------------------------------------	-----

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.004
C	0.015
D	0.059
E	0.234
F	0.938
G	3.75
H	15.0

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature] 16/Jan/2018
Signature/Date